

**Extension Memo**

Product Description: V-PLEX and V-PLEX Plus Human IL-17A Kit
Lot Number: K0082602
New Expiration Date: 30 JUN 2027
Catalog Number(s): K151RF-Series, K156RF-Series

Based on additional testing, the original date of expiration for this kit has been extended by 9 months.

The expiration dates for the following components have also been extended:

Description	Lot Number	Original Expiration Date	New Expiration Date
Human IL-17A 96-Well Small Spot SECTOR Plate	Z0048048	31 MAY 2027	30 JUN 2027
Anti-hu IL-17A Antibody	D0082046	30 SEP 2026	30 JUN 2027
Diluent 3	M0030132	31 MAR 2027	30 JUN 2027
Diluent 43	M0430056	30 SEP 2026	30 JUN 2027

Function	Name	Signature	Date
Quality	Karen Hamilla		25 Aug 2026



CERTIFICATE OF ANALYSIS

Product Description: V-PLEX and V-PLEX Plus Human IL-17A Kit
Lot Number: K0082602
Expiration Date: 30 SEP 2026
Catalog Number: K151RF-Series, K156RF-Series

Lot-Specific Components:

Description	Lot Number	Storage Temperature	Expiration Date
Anti-hu IL-17A Antibody	D0082046	2-8°C	30 SEP 2026
Cytokine Panel 1 (human) Calibrator Blend	A0080383	2-8°C	31 DEC 2028
Human IL-17A 96-Well Small Spot SECTOR Plate	Z0048048	2-8°C	31 MAY 2027
Diluent 3	M0030132	≤ -10°C	31 MAR 2027
Diluent 43	M0430056	≤ -10°C	30 SEP 2026

Additional components may be provided to execute the assay. Additional components are not lot-specific.

Precision, Accuracy, and Sensitivity Specifications:

Spot	Name	Calibrator		Cytokine Panel 1 (human) Control 1		Cytokine Panel 1 (human) Control 2		Cytokine Panel 1 (human) Control 3	
		Recovery	CV	Recovery	CV	Recovery	CV	Recovery	CV
1	IL-17A	80% - 120%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%	90% - 130%	< 20%

Spot	Name	LLOQ			ULOQ		
		Value	Recovery	CV	Value	Recovery	CV
1	IL-17A	3.19	80% - 120%	< 20%	3650	80% - 120%	< 20%

Stock Concentration of Calibrator:

Spot	Name	Conc.	Units
1	IL-17A	6490	pg/mL

Dilution to top of Curve: 1X

This product is intended for Research Use Only. Not for use in Diagnostic Procedures.

	Name	Function	Signature	Date
Review/Approval	Karen Hamilla	Quality		20 FEB 2025